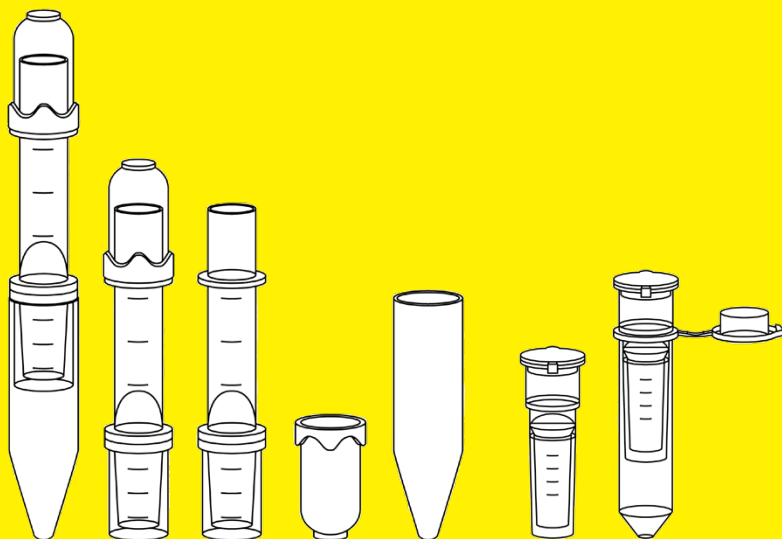


Instructions for Use

Vivaspin[®] 500 and 2

Centrifugal Ultrafiltration Units for General Laboratory Use



3104647-001-00



SARTORIUS

Contents

1	About these Instructions	5
1.1	Scope	5
1.2	Target Groups	7
1.3	Symbols Used	7
1.3.1	Warnings in Operation Descriptions	7
1.3.2	Other Symbols	7
2	Safety Instructions	8
2.1	General Functions	8
2.2	Personnel Qualification	9
2.3	Significance of these Instructions	9
2.4	Functionality of the Product	9
3	Product Description	10
3.1	Vivaspin® 500	10
3.2	Vivaspin® 2	11
3.3	Product Symbols	12
4	Process Preparation	13
4.1	Scope of Delivery	13
4.2	Unpacking	13
5	Operation	14
5.1	Pre-Rinsing the Product	14
5.2	Sanitizing the Product	14
5.3	Applying the Sample	15
5.4	Inserting the Product into the Centrifuge	16
5.5	Performing Filtration	16
5.6	Removing the Sample	16
5.7	Vivaspin® 2	17
5.7.1	Removing the Concentrator from the Filtrate Container	17
5.7.2	Reverse spin with Vivaspin® 2	18
5.8	Desalting or Buffer Exchange	18
6	Storage	19
6.1	Storing the Product	19

7	Disposal	20
7.1	Decontaminating the Product	20
7.2	Disposing of the Product.....	20
8	Technical Specifications	21
8.1	Dimensions	21
8.2	Materials	21
8.3	Ambient Conditions.....	22
8.4	Operating conditions	22
8.4.1	Filtration Volumes	22
8.4.2	Centrifugation Limit Values.....	23
8.5	Equipment Required.....	23
8.5.1	Pipettes.....	23
8.5.2	Centrifuges.....	24
8.6	Sanitizing Methods	24
9	Typical Performance Characteristics	25
9.1	Vivaspin® 500.....	25
9.2	Vivaspin® 2	26
10	Chemical Compatibility	27

1 About these Instructions

1.1 Scope

These instructions are part of the product. These instructions apply to the following versions of the product:

Vivaspin® 500 PES	Quantity	Prod. no.
3 kDa	25 100	VS0191 VS0192
5 kDa	25 100	VS0111 VS0112
10 kDa	25 100	VS0101 VS0102
30 kDa	25 100	VS0121 VS0122
50 kDa	25 100	VS0131 VS0132
100 kDa	25 100	VS0141 VS0142
300 kDa	25 100	VS0151 VS0152
1,000 kDa	25 100	VS0161 VS0162
0.2 µm	25 100	VS0171 VS0172

Vivaspin® 2 PES	Quantity	Prod. no.
3 kDa	25 100	VS0291 VS0292
5 kDa	25 100	VS0211 VS0212
10 kDa	25 100	VS0201 VS0202
30 kDa	25 100	VS0221 VS0222
50 kDa	25 100	VS0231 VS0232
100 kDa	25 100	VS0241 VS0242
300 kDa	25 100	VS0251 VS0252
1,000 kDa	25 100	VS0261 VS0262
0.2 µm	25 100	VS0271 VS0272

Vivaspin® 2 Cellulose Triacetate	Quantity	Prod. no.
10 kDa	25 100	VS02V1 VS02V2
20 kDa	25 100	VS02X1 VS02X2

Vivaspin® 2 Hydrosart®	Quantity	Prod. no.
2 kDa	25 100	VS02H91 VS02H92
5 kDa	25 100	VS02H11 VS02H12
10 kDa	25 100	VS02H01 VS02H02
30 kDa	25 100	VS02H21 VS02H22

1.2 Target Groups

The instructions are designed for the following target groups. The target groups must possess the knowledge listed below.

Target Group	Knowledge and Qualifications
Operator	The operator of the product is responsible for compliance with safety requirements and workplace safety regulations. The operator must ensure that anyone working with the product has access to the relevant information and is trained to work with the product.

1.3 Symbols Used

1.3.1 Warnings in Operation Descriptions

NOTICE

Denotes a hazard that may result in property damage if it is **not** avoided.

1.3.2 Other Symbols

- ▶ Required action: Describes actions that must be carried out. The actions in the sequence must be carried out in succession.
- ▷ Result: Describes the result of the actions carried out.

2 Safety Instructions

2.1 General Functions

The product is intended for the ultrafiltration and | or diafiltration of biological and aqueous solutions. The sample solutions and volumes used must be suitable for the product.

The filtration process must be carried out in a centrifuge. Macromolecules that are sufficiently larger than the nominal pore size of the membrane are retained above the membrane and progressively concentrated. The vertical membrane inhibits membrane fouling while the built-in dead stop impedes concentration to dryness and loss of sample.

The product is supplied non-sterile. It is intended for single use and must be disposed of after one use.

The product is intended exclusively for use in accordance with these instructions. Any further use beyond this is considered improper.

Operating Conditions for the Product

The product is intended for general laboratory use.

The product may only be used with the equipment and under the operating conditions described in the Technical Data section of these instructions.

2.2 Personnel Qualification

Persons without sufficient knowledge in the safe use of the device can injure themselves and others.

If a specific qualification is required for an activity: The target group is indicated. If no qualification is specified: The activity can be carried out by the target group "Operator".

2.3 Significance of these Instructions

Failure to follow the instructions might have serious consequences, e.g. danger to individuals.

- ▶ Read the instructions carefully and completely. The instructions for action build on each other.
- ▶ Ensure that the information contained in these instructions is available to all individuals working with the product.

2.4 Functionality of the Product

A damaged product or worn parts can lead to malfunctions or cause hazards which are difficult to identify.

- ▶ Only operate the product when it is safe and in perfect working order.

3 Product Description

3.1 Vivaspin® 500

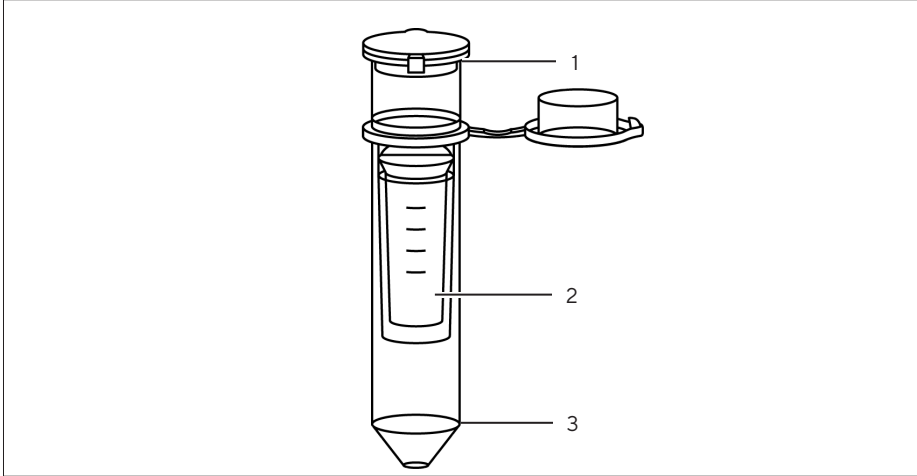


Fig. 1: Product overview (example)

Pos.	Description
1	Concentrator
2	Membrane
3	Filtrate container

3.2 Vivaspin® 2

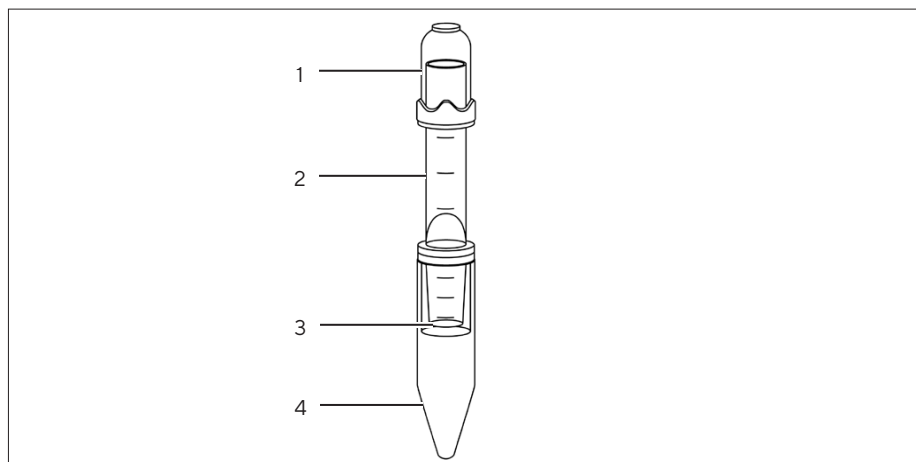









Fig.2: Product overview (example)

Pos.	Description
1	Concentrator cap
2	Concentrator
3	Membrane
4	Filtrate container

3.3 Product Symbols

Symbol	Definition
	Catalogue number
	Do not reuse
	Use by
	Batch code
	Manufacturer
	Temperature limitation
	Non-sterile product

4 Process Preparation

4.1 Scope of Delivery

Article	Qty
Product packed in a cardboard box	
Vivaspin® 500	25 or 100
Vivaspin® 2	25 or 100
Instructions for Use	1

4.2 Unpacking

Procedure

- ▶ **NOTICE** Risk of product malfunctions due to exceeding the usability!
Check the usability of the product (see specification on packaging).
Dispose of products for which the usability has been exceeded.
- ▶ Unpack the product.

5 Operation

5.1 Pre-Rinsing the Product

Membranes in the product may contain traces of glycerin. If this substance can interfere with the analysis of the sample: The membrane may be rinsed before filtration.

Procedure

- ▶ Remove the concentrator cap.
- ▶ Use a pipette to apply a filling volume of buffer solution or deionized water into the concentrator.
- ▶ Replace the concentrator cap.
- ▶ Wash the buffer solution or deionized water through the membrane by centrifugation.
- ▶ Empty the concentrator and filtrate container.
- ▶ If the pre-rinsed product is not used immediately: Cover the surface of the membrane with buffer solution or water and store the product in the refrigerator. The membrane must not dry out.

5.2 Sanitizing the Product

The product can be sanitized before use. The sanitizing method must be suitable for the product (see Chapter "8.6 Sanitizing Methods," page 24).

Procedure

- ▶ Remove the concentrator cap.
- ▶ Sanitize the product using the desired sanitizing method.
- ▶ Empty the product.

5.3 Applying the Sample

It is recommended that a pipette is used to apply the sample into the product. The pipette must be compatible with the product (see Chapter "8.5.1 Pipettes," page 23).

Please ensure that the molecular weight cut-off (MWCO) of the product is suitable for the size of the target molecule to be concentrated. In order to ensure maximum recovery of the target molecule, it is recommended to select a MWCO that is at least 50% below the size of the target molecule.

NOTICE

Risk of product malfunctions due to using unsuitable samples!

- ▶ Only pour suitable samples into the product (see Chapter "10 Chemical Compatibility," page 27).

NOTICE

Risk of product malfunctions or damage to the centrifuge due to exceeding the maximum filling volume!

- ▶ Do **not** exceed the maximum filling volume (see Chapter "8.4.1 Filtration Volumes," page 22).

Procedure

- ▶ Check whether the MWCO of the product is suitable for the application.
- ▶ Remove the concentrator cap.
- ▶ Apply the sample into the product using a pipette. Comply with the maximum filling volume.
- ▶ Replace the concentrator cap.

5.4 Inserting the Product into the Centrifuge

Procedure

- ▶ **NOTICE** Risk of product malfunctions or damage to the centrifuge! Only use the product in suitable centrifuges (see Chapter "8.5.2 Centrifuges," page 24).
- ▶ Insert the product into the centrifuge.
- ▶ If a centrifuge with fixed-angle rotor is used: Place the product into the centrifuge so that the printed volume graduations on the concentrator are facing upwards.

5.5 Performing Filtration

Procedure

- ▶ **NOTICE** Risk of product malfunctions or damage to the centrifuge. Comply with the approved centrifugation limit values (see Chapter "8.4.2 Centrifugation Limit Values," page 23).
- ▶ Centrifuge the product in the centrifuge until the desired concentration level is achieved.

5.6 Removing the Sample

Procedure

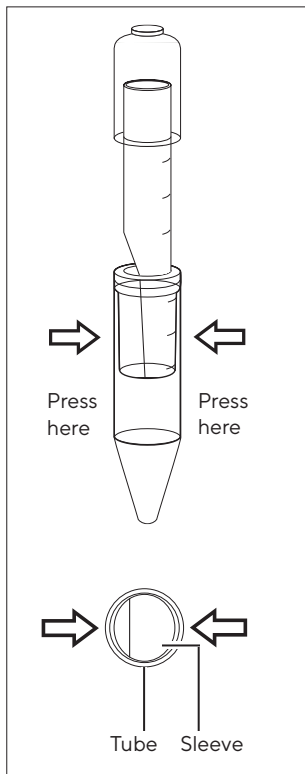
- ▶ If the filtration or concentration is complete: Remove the product from the centrifuge.
- ▶ Remove the concentrator cap.
- ▶ Recover the sample from dead stop pocket of the concentrator using a pipette.
- ▶ If the membrane was pre-rinsed before filtration: Decant the filtrate and concentrate.

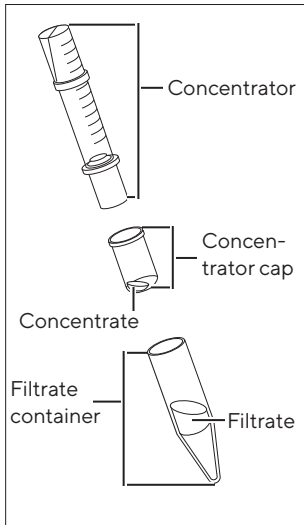
5.7 Vivaspin® 2

5.7.1 Removing the Concentrator from the Filtrate Container

Procedure

- ▶ To release the tube from the concentrator, pinch the tube - to press it into an oval shape - before removing it with a twisting action.





5.7.2 Reverse spin with Vivaspin® 2

The concentrate can be collected into the concentrator cap by reverse spinning, instead of using a pipette.

Procedure

- ▶ Remove filtrate container.
- ▶ With the concentrator cap still fitted, invert the concentrator.
- ▶ Insert concentrator cap into filtrate container.
- ▶ Insert the product into the centrifuge.
- ▶ If a centrifuge with fixed-angle rotor is used: Place the product into the centrifuge so that the printed volume graduations on the concentrator are facing upwards.
- ▶ Spin at up to 3,000 g for 2 minutes.

5.8 Desalting or Buffer Exchange

Procedure

- ▶ Concentrate the sample to the desired level.
- ▶ Remove the concentrator cap.
- ▶ Discard the filtrate.
- ▶ Refill the concentrator with an appropriate exchange buffer.
- ▶ Concentrate the sample again.
- ▶ Repeat the process until the original buffer and | or contaminating microsolite has been sufficiently removed.
- ▶ If the desalting or buffer exchange is complete: Recover the sample.

6 Storage

6.1 Storing the Product

If the product has been unpacked and membrane has been pre-rinsed: The membrane must be protected against drying out. For this purpose, the membrane must be stored in a moist and cool condition.

NOTICE

Risk of damage to the product due to improper storage!

- ▶ Comply with the storage specifications.
-

Procedure

- ▶ If the product is packaged: Store the product in the packaging.
- ▶ If the product has been unpacked and the membrane has been pre-rinsed:
 - ▶ Remove the concentrator cap.
 - ▶ Cover the membrane with buffer solution or water.
 - ▶ Replace the concentrator cap.
- ▶ Store the product according to the ambient conditions (see Chapter "8.3 Ambient Conditions," page 22).

7 Disposal

7.1 Decontaminating the Product

If the product has come into contact with hazardous substances: Steps must be taken to ensure proper decontamination and declaration. The operator of the product is responsible for adhering to local government regulations on the proper decontamination and declaration for transport and disposal.

Procedure

- ▶ If the product has come into contact with hazardous substances: Decontaminate the product.

7.2 Disposing of the Product

The product must be disposed of properly. The packaging is made of environmentally friendly materials that can be used as secondary raw materials.

Requirements

The product must be decontaminated.

Procedure

- ▶ Dispose of the product in accordance with local government regulations.
- ▶ Dispose of the packaging in accordance with local government regulations.

8 Technical Specifications

8.1 Dimensions

	Vivaspin® 500		Vivaspin® 2	
	Unit	Value	Unit	Value
Length x Diameter	mm	50 x 11	mm	126 x 17
Active membrane surface	cm ²	0.5	cm ²	1.2
Weight	g	3	g	11

8.2 Materials

	Vivaspin® 500		Vivaspin® 2	
	Unit	Value	Unit	Value
Concentrator		Polycarbonate		Polycarbonate
Filtrate container		Polypropylene		Polycarbonate
Membrane		Polyethersulfone		Polyethersulfone, Cellulose triacetate, Hydrosart®

8.3 Ambient Conditions

	Unit	Value
Storage temperature		
When packed	°C	+15 - +30
When unpacked, with membrane kept moist	°C	+2 - +8

8.4 Operating conditions

8.4.1 Filtration Volumes

	Unit	Centrifuge with swing bucket rotor	Centrifuge with fixed-angle rotor (25°)
	Unit	Value	Value
Vivaspin® 500			
Filling volume, maximum	µL	-	500
Membrane hold-up volume, minimum	µL	-	< 5
Dead stop volume ¹	µL	-	5
Vivaspin® 2			
Filling volume, maximum	mL	3	2
Membrane hold-up volume, minimum	µL	< 10	< 10
Dead stop volume ¹	µL	8	8

¹The dead stop volume may vary depending on the type and concentration of the sample, operating temperature and | or centrifuge rotor

8.4.2 Centrifugation Limit Values

		Centrifuge with swing bucket rotor	Centrifuge with fixed angle rotor
	Unit	Value	Value
Vivaspin® 500	<i>g</i>	-	12,000
Vivaspin® 2	<i>g</i>	4,000	8,000

8.5 Equipment Required

8.5.1 Pipettes

Pasteur pipette, variable volume or fixed volume pipette for sample application and concentrate or filtrate retrieval.

8.5.2 Centrifuges

Centrifuge accepting conical base tubes.

	Required carriers	
	Unit	Value
Vivaspin® 500		
Rotor type	Fixed angle	
Volume	mL	2.2
Diameter	mm	11
Minimum rotor angle	°	40
Vivaspin® 2		
Rotor type	Swing bucket or fixed angle	
Volume	mL	15
Diameter	mm	17
Minimum rotor angle	°	25

8.6 Sanitizing Methods

Rinsing with 70% ethanol or with sanitizing gas mixture, e.g. ethylene oxide

Not suitable for autoclaving

9 Typical Performance Characteristics

9.1 Vivaspin® 500

	Time to concentrate up to 30x at 20°C	
Start volume	500 µL	
	Time (min)	Solute Recovery
Aprotinin 0.25 mg/mL (6.5 kDa)		
3 kDa PES	30	96%
BSA 1.0 mg/mL (66 kDa)		
5 kDa PES	15	96%
10 kDa PES	5	96%
30 kDa PES	5	95%
IgG 0.25 mg/mL (160 kDa)		
30 kDa PES	10	96%
50 kDa PES	10	96%
100 kDa PES	10	96%

9.2 Vivaspin® 2

Time to concentrate up to 30x at 20°C		
Start volume	2 mL	
	Time (min)	Solute Recovery
Insulin chain A 0.1 mg/mL (2.5 kDa)		
2 kDa Hydrosart®	35	95%
Aprotinin 0.25 mg/mL (6.5 kDa)		
3 kDa PES	50	96%
BSA 1.0 mg/mL (66 kDa)		
5 kDa PES	12	98%
5 kDa Hydrosart®	22	98%
10 kDa PES	8	98%
10 kDa CTA	10	96%
10 kDa Hydrosart®	12	98%
20 kDa CTA	5	96%
30 kDa PES	8	97%
30 kDa Hydrosart®	5	97%
IgG 0.25 mg/mL (160 kDa)		
20 kDa CTA	6	97%
30 kDa PES	10	96%
50 kDa PES	10	96%
100 kDa PES	8	95%

10 Chemical Compatibility

Chemical, biological and aqueous solutions with appropriate compatibility for the materials of the product (2 hr contact time)

Solutions	PES	CTA	HY
Compatible pH range	pH 1-9	pH 4-8	pH 1-9
Acetic Acid (25%)	OK	NO	OK
Acetone (10%)	NO	NO	NO
Acetonitrile (10%)	NO	NO	NO
Ammonium Hydroxide (5%)	?	OK	OK
Ammonium Sulphate (saturated)	OK	?	?
Benzene (100%)	NO	NO	NO
n-Butanol (70%)	?	NO	?
Chloroform (1%)	NO	NO	NO
Dimethyl Formamide (10%)	?	NO	NO
Dimethyl Sulfoxide (5%)	OK	NO	NO
Ethanol (70%)	OK	OK	OK
Ethyl Acetate (100%)	NO	NO	NO
Formaldehyde (30%)	OK	OK	OK
Formic Acid (5%)	OK	?	OK
Glycerine (70%)	OK	OK	OK
Guanidine HCl (6 M)	OK	?	OK
Hydrocarbons, aromatic	NO	NO	NO
Hydrocarbons, chlorinated	NO	NO	NO
Hydrochloric Acid (1 M)	OK	NO	OK

Chemical, biological and aqueous solutions with appropriate compatibility for the materials of the product (2 hr contact time)

Solutions	PES	CTA	HY
Compatible pH range	pH 1-9	pH 4-8	pH 1-9
Imidazole (300 mM)	OK	NO	?
Isopropanol (70%)	OK	OK	OK
Lactic Acid (5%)	OK	NO	OK
Mercaptoethanol (1 M)	NO	NO	OK
Methanol (60%)	?	?	OK
Nitric Acid (10%)	OK	NO	NO
Phenol (1%)	?	?	NO
Phosphate Buffer (1 M)	OK	OK	OK
Polyethylene Glycol (10%)	OK	?	?
Pyridine (100%)	NO	NO	NO
Sodium Carbonate (20%)	OK	NO	?
Sodium Deoxycholate (5%)	OK	?	?
Sodium Dodecylsulfate (0.1 M)	OK	OK	OK
Sodium Hydroxide (2.5 M)	NO	NO	NO
Sodium Hypochlorite (200 ppm)	OK	NO	NO
Sodium Nitrate (1%)	OK	?	OK
Sulfamic Acid (5%)	OK	NO	?
Tetrahydrofuran (5%)	NO	NO	NO
Toluene (1%)	NO	NO	NO
Trifluoroacetic Acid (10%)	OK	NO	OK
Tween®* 20 (0.1%)	OK	OK	OK

Chemical, biological and aqueous solutions with appropriate compatibility for the materials of the product (2 hr contact time)

Solutions	PES	CTA	HY
Compatible pH range	pH 1-9	pH 4-8	pH 1-9
Triton ^{®**} X-100 (0.1%)	OK	OK	OK
Urea (8 M)	OK	?	OK

OK = Acceptable ? = Questionable NO = Not recommended

* Triton[®] is a registered trademark of Union Carbide Corp.

** Tween[®] is a registered trademark of ICI Americas Inc.

Sartorius Stedim Lab Ltd.
Sperry Way, Stonehouse
GL10 3UT, UK

Phone: +44 1453 821972
www.sartorius.com

The information and figures contained in these instructions correspond to the version date specified below.

Sartorius reserves the right to make changes to the technology, features, specifications and design of the equipment without notice.

Masculine or feminine forms are used to facilitate legibility in these instructions and always simultaneously denote all genders.

Copyright notice:

These instructions, including all components, are protected by copyright.

Any use beyond the limits of the copyright law is not permitted without our approval.

This applies in particular to reprinting, translation and editing irrespective of the type of media used.

Last updated:

05 | 2023

© 2023
Sartorius Stedim Lab Ltd.
Sperry Way, Stonehouse
GL10 3UT, UK

KS | Publication No.: SLU6093-e230513